

PRODUCT INFORMATION



Lythrine

Item No. 25804

CAS Registry No.: 5286-10-2
Formal Name: (4a*S*,6*S*,9*Z*,20*S*)-2,3,4,4a,5,6-hexahydro-14-hydroxy-17,18-dimethoxy-20H-6,20-methano-11,15-metheno-1*H*,8*H*-benzo[*g*]pyrido[2,1-*d*][1,5]oxaazacyclohexadecin-8-one

Synonym: (+)-Lythrine

MF: C₂₆H₂₉NO₅

FW: 435.5

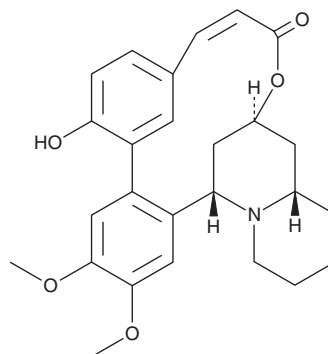
Purity: ≥95%

UV/Vis.: λ_{max}: 284 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Lythrine is supplied as a crystalline solid. A stock solution may be made by dissolving the lythrine in the solvent of choice. Lythrine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of lythrine in ethanol is approximately 1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

Lythrine is an alkaloid originally isolated from *H. salicifolia* that inhibits prostaglandin synthetase (IC₅₀ = 469 μM).¹

Reference

1. Lema, W.J., Blankenship, J.W., and Malone, M.H. Prostaglandin synthetase inhibition by alkaloids of *Heimia salicifolia*. *J. Ethnopharmacol.* **15(2)**, 161-167 (1986).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/17/2018

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM